

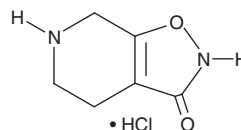
PRODUCT INFORMATION



Gaboxadol (hydrochloride)

Item No. 16355

CAS Registry No.: 85118-33-8
Formal Name: 4,5,6,7-tetrahydro-isoxazolo[5,4-c]pyridin-3(2H)-one, monohydrochloride
Synonym: THIP
MF: C₆H₈N₂O₂ • HCl
FW: 176.6
Purity: ≥98%
UV/Vis.: λ_{max}: 210 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Gaboxadol (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the gaboxadol (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Gaboxadol (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of gaboxadol (hydrochloride) in these solvents is approximately 20 and 0.1 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of gaboxadol (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of gaboxadol (hydrochloride) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Gaboxadol is a GABA_A receptor agonist whose potency varies depending on the receptor subunit composition (partial agonist at α1β2γ2 (ED₅₀ = 143 μM), full agonist at α5 (ED₅₀ = 28-129 μM), and super agonist at α4β3δ (ED₅₀ = 6 μM)).¹ It also acts as an antagonist at ρ1 GABA_C receptors (IC₅₀ = 25 μM).¹ Gaboxadol has been studied as a non-opioid analgesic and a novel hypnotic agent.²⁻⁴

References

1. Johnston, G.A.R., Chebib, M., Hanrahan, J.R., *et al.* Neurochemicals for the investigation of GABA_C receptors. *Neurochem. Res.* **35(12)**, 1970-1977 (2010).
2. Drasbek, K.R. and Jensen, K. THIP, a hypnotic and antinociceptive drug, enhances an extrasynaptic GABA_A receptor-mediated conductance in mouse neocortex. *Cereb. Cortex* **16(8)**, 1134-1141 (2006).
3. Vashchinkina, E., Panhelainen, A., Vekovischeva, O.Y., *et al.* GABA site agonist gaboxadol induces addiction-predicting persistent changes in ventral tegmental area dopamine neurons but is not rewarding in mice or baboons. *J. Neurosci.* **32(15)**, 5310-5320 (2012).
4. Deacon, S., Staner, L., Staner, C., *et al.* Effect of short-term treatment with gaboxadol on sleep maintenance and initiation in patients with primary insomnia. *Sleep* **30(3)**, 281-287 (2007).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/15/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM